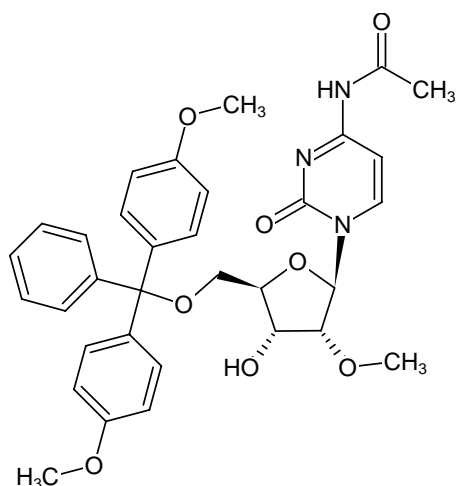


**5'-DMT-2'-(O-methyl)-cytidine (N-Ac)**

Cat. No.	Amount
N-PM-5735-500	500 mg
N-PM-5735-1G	1 g



Structural formula of 5'-DMT-2'-(O-methyl)-cytidine (N-Ac)

For research use only!**Shipping:** shipped at ambient temperature**Storage Conditions:** store at -20 °C**Shelf Life:** 30 months after date of delivery**Molecular Formula:** C₃₃H₃₅N₃O₈**Molecular Weight:** 601.65 g/mol**Purity:** ≥ 98 % (HPLC) (Isomeric impurity 0 %)**Form:** white crystalline powder**Solubility:** acetonitrile